

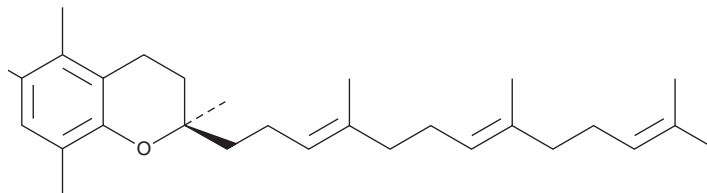
PRODUCT INFORMATION



β -Tocotrienol

Item No. 9001533

CAS Registry No.: 490-23-3
Formal Name: (2R)-3,4-dihydro-2,5,8-trimethyl-2-[4,8,12-trimethyl-3E,7E,11-tridecatrien-1-yl]-2H-1-benzopyran-6-ol
MF: C₂₈H₄₂O₂
FW: 410.6
Purity: \geq 98%
UV/Vis.: λ_{max} : 293 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: \geq 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

β -Tocotrienol is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of β -tocotrienol in these solvents is approximately 10 mg/ml.

β -Tocotrienol is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

The four tocopherols (α , β , δ , and γ) and four tocotrienols (α , β , δ , and γ) are forms of vitamin E, which is known for its antioxidant activity.¹ Tocotrienols are found in high concentrations in palm oil and wheat germ. β -Tocotrienol is a less potent antioxidant than α -tocotrienol (Item No. 10008377).¹ A high cholesterol diet supplemented with a mixture of α - and β -tocopherols and tocotrienols has been shown to lower hepatic cholesterol and cholesterol ester concentration and to reduce atherosclerotic lesions in ApoE^{+/-} mice.²

References

1. Kamal-Eldin, A. and Appelqvist, L.-Å. The chemistry and antioxidant properties of tocopherols and tocotrienols. *Lipids* **31**, 671-701 (1996).
2. Black, T.M., Wang, O., Maeda, N., et al. Palm tocotrienols protect ApoE +/- mice from diet-induced atheroma formation. *J. Nutr.* **130(10)**, 2420-2426 (2000).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 09/28/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM